

MEQ [6-Methoxy-N-ethylquinolinium iodide]Catalog number: 21250
Unit size: 100 mg**Product Details**

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

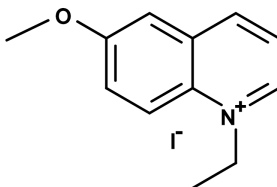
Chemical Properties

Appearance White solid

Molecular Weight 315.15

Soluble In DMSO

Chemical Structure

**Spectral Properties**

Excitation Wavelength 318 nm

Emission Wavelength 447 nm

Applications

A chloride indicator that has the fluorescence dependent of chloride concentration. Its fluorescence is quenched by chloride through collision quenching.